

| | |
|-----------------|---------------|
| Product Number: | 102153 |
| Lot Number: | 501797 |
| Expiration: | 1-Apr-2020 |
| Platform(s): | HD-1 and SR-X |



Data below represents results generated on the Simoa™ HD-1 Analyzer.

| Component | Part Number | Lot Number |
|----------------------|---------------|------------------|
| Bead Reagent | 102141 | 912607 |
| Detector Reagent | 102143 | 912606 |
| SBG Reagent | 102145 | 911408 |
| Sample Diluent | 102147 | 911407 |
| Calibrators | 102150 | 905901 |
| Controls | 102151-102152 | 905902 |
| SBG Diluent (SR-X) † | 103158 | N/A ¹ |
| RGP Reagent | 103159 | N/A ¹ |

¹ Reagents are not Kit Lot Specific

† Specific for use on the SR-X platform - refer to the Simoa™ Neuro 4-Plex A SR-X Kit Instructions (KI-0031).

Release Materials

Tau

| | Lot | Result (pg/mL) | Mean Range (pg/mL) |
|-----------|--------------|----------------|--------------------|
| Control 1 | 905902 | 2.07 | 1.59-2.38 |
| Control 2 | 905902 | 93.9 | 76.2-114 |
| Panel 1 | MF-101617-01 | 1.68 | 0.650-1.79 |
| Panel 2 | MF-101617-02 | 36.0 | 22.7-44.9 |
| Panel 3 | MF-101617-03 | 140 | 93.3-204 |

NF-light®

| | Lot | Result (pg/mL) | Mean Range (pg/mL) |
|-----------|--------------|----------------|--------------------|
| Control 1 | 905902 | 5.64 | 3.90-6.20 |
| Control 2 | 905902 | 562 | 401-602 |
| Panel 1 | MF-101617-01 | 5.22 | 3.08-6.35 |
| Panel 2 | MF-101617-02 | 41.5 | 16.8-50.2 |
| Panel 3 | MF-101617-03 | 400 | 232-548 |

GFAP®

| | Lot | Result (pg/mL) | Mean Range (pg/mL) |
|-----------|--------------|----------------|--------------------|
| Control 1 | 905902 | 23.3 | 18.0-27.0 |
| Control 2 | 905902 | 1242 | 912-1369 |
| Panel 1 | MF-101617-01 | 49.0 | 40.6-62.2 |
| Panel 2 | MF-101617-02 | 181 | 110-229 |
| Panel 3 | MF-101617-03 | 2332 | 2138-3313 |

UCHL1®

| | Lot | Result (pg/mL) | Mean Range (pg/mL) |
|-----------|--------------|----------------|--------------------|
| Control 1 | 905902 | 119 | 65.7-136 |
| Control 2 | 905902 | 10311 | 7350-11026 |
| Panel 1 | MF-101617-01 | 70.6 | 39.0-89.1 |
| Panel 2 | MF-101617-02 | 318 | 308-473 |
| Panel 3 | MF-101617-03 | 4655 | 4253-6380 |

Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Review/Approval

Patrizia Stadler QC Assoc. Director

Name

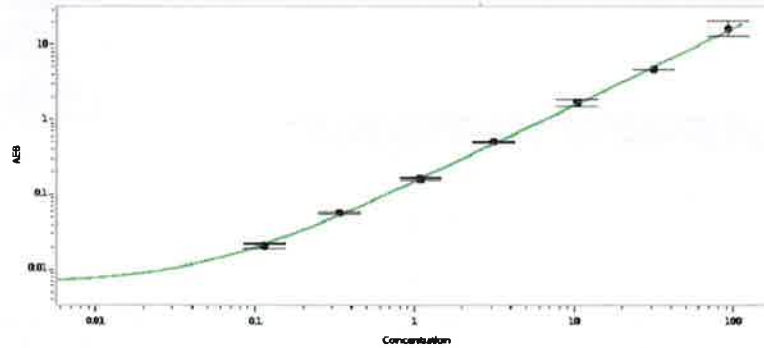
Title

P. Stadler

09 MAY 2019

Signature/Date

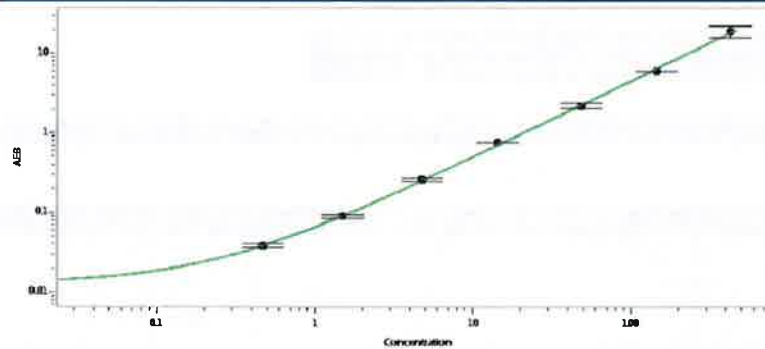
Tau Calibration Curve



Calibrator Levels (pg/mL)

| | |
|---|-------|
| A | 0.000 |
| B | 0.113 |
| C | 0.332 |
| D | 1.08 |
| E | 3.12 |
| F | 10.3 |
| G | 31.3 |
| H | 91.6 |

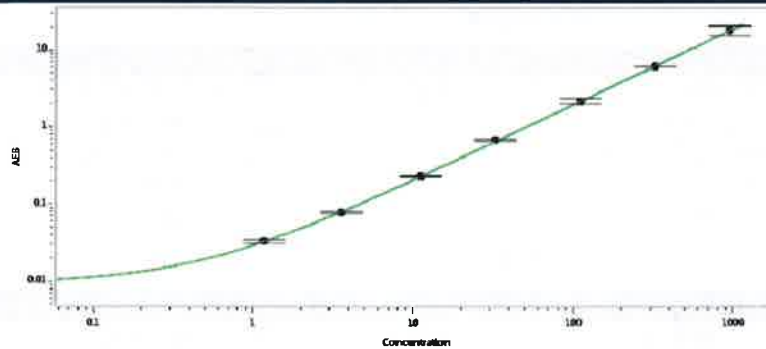
NF-light® Calibration Curve



Calibrator Levels (pg/mL)

| | |
|---|-------|
| A | 0.000 |
| B | 0.470 |
| C | 1.50 |
| D | 4.80 |
| E | 14.3 |
| F | 48.0 |
| G | 144 |
| H | 426 |

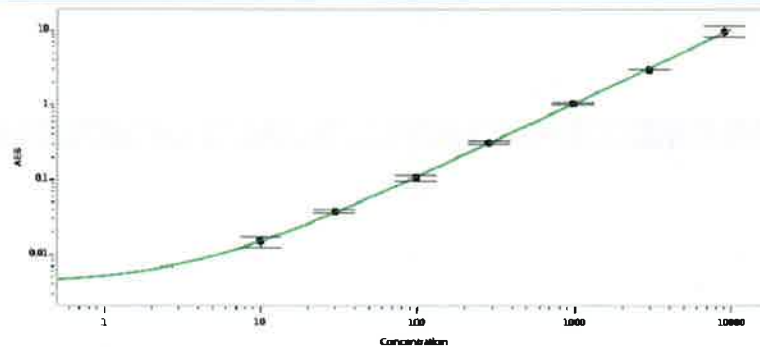
GFAP® Calibration Curve



Calibrator Levels (pg/mL)

| | |
|---|-------|
| A | 0.000 |
| B | 1.16 |
| C | 3.54 |
| D | 11.1 |
| E | 32.6 |
| F | 111 |
| G | 323 |
| H | 946 |

UCHL1® Calibration Curve



Calibrator Levels (pg/mL)

| | |
|---|-------|
| A | 0.000 |
| B | 10.0 |
| C | 29.6 |
| D | 97.9 |
| E | 285 |
| F | 980 |
| G | 2940 |
| H | 8998 |